

n-Octyl-β-D-glucopyranoside (analytical grade)

Non-ionic detergent. Highly purified.

Highly purified nonionic detergent that solubilizes membrane-bound proteins in their native state.

Citations: 5

[View Online »](#)

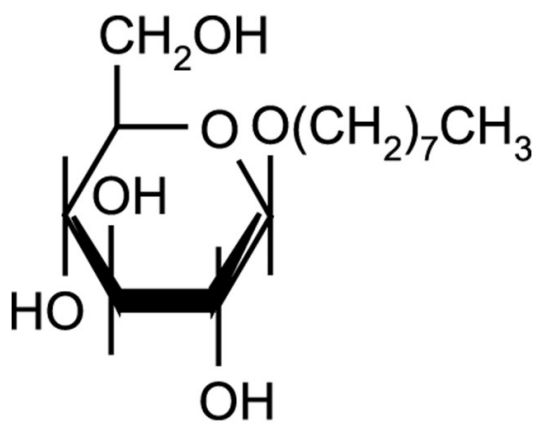
Ordering Information

[Order Online »](#)

ALX-500-001-G001	1g
ALX-500-001-G005	5g
ALX-500-001-G025	25g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep under inert gas. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	OG
Appearance	White to off-white powder.
CAS	29836-26-8
Formula	$C_{14}H_{28}O_6$
Identity	Determined by 1H -NMR.
MW	292.4
Purity	≥99% (HPLC)
Solubility	Soluble in water (100mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)